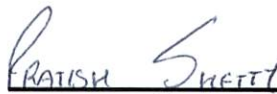


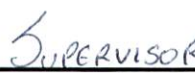
Product Number:	103529
Lot Number:	231108
Expiration:	7-Nov-2023
Storage:	2 - 8°C


Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.38 x 10 ⁹ Beads/mL

Review/Approval


 Name

QC

 SUPERVISOR
 Title


 Signature/Date

20 APR 2023